

# Complement component C1s (human) monoclonal antibody (9)

C1s proenzyme is a 79.8kDa glycoprotein produced primarily by hepatocytes but also by other cells. The proenzyme is produced as a single-chain polypeptide of 688 amino acids. Upon activation, the peptide is cleaved between R437 and I438, producing an activated enzyme consisting of two disulfide-linked chains, A chain (52.2kDa) and B chain (27.7kDa). The calcium-dependent catalytic subunit of the C1 complex (C1q, C1r, C1s) is comprised of two C1s and two C1r molecules, and this serine protease tetramer is responsible for the initial steps in the cleavage of C4 and C2 in the classical complement pathway.

This antibody is covered by our [Worry-Free Guarantee](#).

## Ordering Information

[Order Online »](#)

BPD-ABS-002-09-1	1mg
BPD-ABS-002-09-04	400µg

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** C1 esterase, Complement C1s subcomponent, C1s proenzyme

**Application** ELISA, IHC

**Clone** 9

**Formulation** Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.

**GenBank ID** 716

**Host** Mouse

**Immunogen** Human plasma complement component C1s Proenzyme.

**Isotype** IgG1κ

**Purity Detail** Protein A/G-purified.

**Source** From culture supernatant.

**Species Reactivity** Human

**Specificity** Specific for C1s proenzyme and activated free C1s.

**UniProt ID** P09871

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)